



# MetaMorph® NX Microscopy Automation and Image Analysis Software

## Complexity made simple!

### Easy-to-Find Tools

- An interface with the user in mind. Find the right tool quickly and complete the task; simplifies training.
- Interactive zoom and pan within an entire image dataset. Easily focus in on the most important results.

### Workflow Oriented

- Consistent workflow-oriented interface guides you through even the most complicated acquisition and analysis tasks.
- Minimizes number of steps required for experiment setup without the need for complex macros or system configuration.
- Acquired datasets are easily analyzed; tabular and graphical extracted information remains linked to image source.

### Clean Desktop

- Patent-pending desktop is always clean and organized for a clear view of your data.
- Clearly visualize and compare multi-parametric datasets across different data views.
- Quickly access views optimized for time, position, and wavelength.

### Experiment History

- Experiment aware, all images, measurements and acquisition parameters stay together throughout experiment cycle.

